

PRODUCT INFORMATION



Pravastatin-d₃ (sodium salt)

Item No. 25464

CAS Registry No.: 1329836-90-9
Formal Name: (3R,5R)-3,5-dihydroxy-7-
((1S,2S,6S,8S,8aR)-6-hydroxy-2-methyl-
8-(((S)-2-(methyl-d₃)butanoyl)oxy)-
1,2,6,7,8,8a-hexahydronaphthalen-1-yl)
heptanoate, monosodium salt

MF: C₂₃H₃₂D₃O₇ • Na
FW: 449.5

Chemical Purity: ≥98% (Pravastatin)

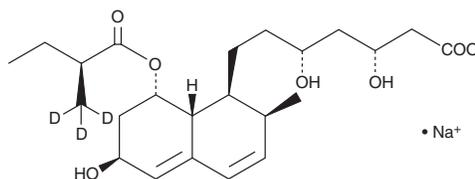
Deuterium

Incorporation: ≥99% deuterated forms (d₁-d₃); ≤1% d₀

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Pravastatin-d₃ is intended for use as an internal standard for the quantification of pravastatin (Item No. 10010343) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Pravastatin-d₃ (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the pravastatin-d₃ (sodium salt) in the solvent of choice, which should be purged with an inert gas. Pravastatin-d₃ (sodium salt) is slightly soluble in methanol.

Description

Pravastatin is a potent and competitive HMG-CoA reductase inhibitor ($K_i = 2.3$ nM for the active, open ring form of the molecule) that is a ring-hydroxylated metabolite of mevastatin (Item No. 10010340).¹ *In vivo*, pravastatin reduces total plasma cholesterol levels by 29% in dogs when administered at a dose of 20 mg/kg per day for five weeks. Formulations containing pravastatin have been used to reduce LDL cholesterol and triglyceride levels and increase HDL cholesterol in the prevention of cardiovascular disease.

Reference

1. Corsini, A., Maggi, F.M., and Catapano, A.L. Pharmacology of competitive inhibitors of HMG-CoA reductase. *Pharmacol. Res.* **31**(1), 9-27 (1995).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 05/04/2023

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM